

PRODUCT INFORMATION

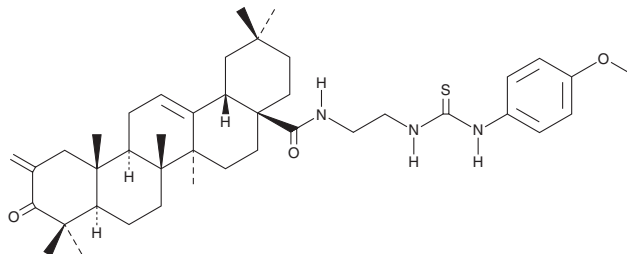


CAY10736

Item No. 26754

CAS Registry No.: 2251753-61-2
Formal Name: N-[2-[[[(4-methoxyphenyl)amino]thioxomethyl]amino]ethyl]-2-methylene-3-oxo-olean-12-en-28-amide

Synonym: Compound 51
MF: C₄₁H₅₉N₃O₃S
FW: 674.0
Purity: ≥90%
UV/Vis.: λ_{max}: 244 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CAY10736 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10736 in the solvent of choice, which should be purged with an inert gas. CAY10736 is soluble in organic solvents such as chloroform and ethanol.

Description

CAY10736 is an anticancer compound.¹ It inhibits proliferation in a panel of melanoma and breast, pancreatic, and lung cancer cell lines (IC₅₀s = 0.827-9.89 μM). CAY10736 inhibits migration and epithelial-to-mesenchymal transition (EMT) of A375 and B16/F10 melanoma cells *in vitro* in a concentration-dependent manner. It reduces the viability of spheroid A375 and B16/F10 cells (IC₅₀s = 2.4 and 1.59 μM, respectively) and increases the production of reactive oxygen species (ROS) in these cells in a concentration-dependent manner. CAY10736 (5 mg/kg) reduces tumor growth in B16/F10 melanoma and Lewis lung carcinoma mouse models and an A375 mouse xenograft model.

Reference

1. Liu, X., Li, B., Zhang, Z., *et al.* Synthesis and discovery novel anti-cancer stem cells compounds derived from the natural triterpenoid acids. *J. Med. Chem.* **61**(23), 10814-10833 (2018).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/16/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM